

DMNPE-Caged D-Luciferin

DMNPE-caged D-luciferin (D-Luciferin, 1-(4,5-dimethoxy-2-nitrophenyl)ethyl ester) is a D-luciferin ester derivative which can cross cell membranes efficiently.^{1,2} Once inside the cells, the ester is continuously hydrolyzed to a supply of D-luciferin.



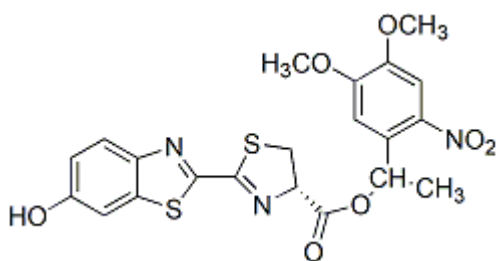
Product attributes

CAS number	223920-67-0
------------	-------------

Product Description

DMNPE-caged D-luciferin (D-Luciferin, 1-(4,5-dimethoxy-2-nitrophenyl)ethyl ester) is a D-luciferin ester derivative which can cross cell membranes efficiently.^{1,2} Once inside the cells, the ester is continuously hydrolyzed to a supply of D-luciferin. Alternatively a burst of D-luciferin is generated by UV photolysis.

- Light yellow solid soluble in DMSO
- Store at -20 °C and protect from light
- C₂₁H₁₉N₃O₇S₂
- MW: 489.52



References

1. Lett Appl Microbiol 30, 223 (2000).
2. Biotechniques 15, 848 (1993).

This datasheet was generated on June 23, 2026 at 06:07:05 PM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/dmnpe-caged-d-luciferin-d-luciferin-1-4-5-dimethoxy-2-nitrophenylethyl-ester/>